

**Purified Anti-Human/monkey
Apolipoprotein E (ApoE) Antibody *E981***Catalog number: V101090
Unit size: 250 µg**Product Details**

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide

Antibody Properties

Species Reactivity	Nhp, human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG2a
Immunogen	apoE (APOE, AD2, APO-E, LDLCQ5, LPG, apolipoprotein E, ApoE4)
Clone	E981

Biological Properties

Preparation	Antibody purified by affinity chromatography
Application	Flow Cytometry (FACS)

Applications
